

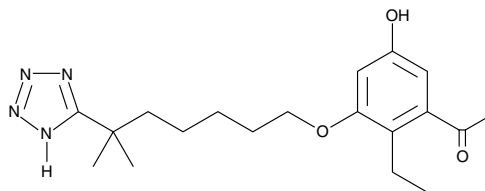
Product Information



LY2552833

Item No. 70715

CAS Registry No.: 117690-79-6
Formal Name: 1-[5-ethyl-2-hydroxy-4-[[6-methyl-6-(1H-tetrazol-5-yl)heptyl]oxy]phenyl]-ethanone
MF: C₁₉H₂₈N₄O₃
FW: 360.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 213, 232, 276, 326 nm



Laboratory Procedures

For long term storage, we suggest that LY255283 be stored as supplied at -20°C. It should be stable for at least one year.

LY255283 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY255283 in an organic solvent purged with an inert gas. LY255283 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of LY255283 in these solvents is approximately 1, 30, and 50 mg/ml, respectively.

LY255283 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY255283 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LY255283 has a solubility of approximately 100 µg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Antagonists of the chemotactic and inflammatory lipoxygenase product leukotriene B₄ (LTB₄) have been potential drug development targets for several years.^{1,2} The tetrazole LY255283 is a competitive antagonist of the BLT₂ receptor. It displaces radiolabeled LTB₄ from guinea pig lung membrane, with an IC₅₀ of about 100 nM.³ LY255283 exhibits IC₅₀ values of ~950 nM and >10 µM at human recombinant BLT₂ and BLT₁ receptors, respectively.⁴ LY255283 inhibits eosinophil chemotaxis by 80% at a concentration of 10 µM, and inhibits the binding of radiolabeled LTB₄ to eosinophil membranes with an IC₅₀ of 260 nM.¹

References

1. Richards, I.M., Sun, F.F., Taylor, B.M., *et al.* Contribution of leukotriene B₄ to airway inflammation and the effect of antagonists. *Ann. N. Y. Acad. Sci.* **629**, 274-287 (1991).
2. Taylor, B.M., Crittenden, N.J., Bruden, M.N., *et al.* Biological activity of leukotriene B₄ analogs: Inhibition of guinea pig eosinophil migration *in vitro* by 2,6-disubstituted pyridine analogs U-75,302 and U-75,485. *Prostaglandins* **42**, 211-224 (1991).
3. Silbaugh, S.A., Stengel, P.W., Cockerham, S.L., *et al.* Pulmonary actions of LY255283, a leukotriene B₄ receptor antagonist. *Eur. J. Pharmacol.* **223**, 57-64 (1992).
4. Yokomizo, T., Kato, K., Hagiya, H., *et al.* Hydroxyicosanoids bind to and activate the low affinity leukotriene B₄ receptor, BLT₂. *J. Biol. Chem.* **276(15)**, 12454-12459 (2001).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/70715

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/19/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com